

HP1 β antibody (mAb)

Catalog Nos: 39979, 39980

RRID: AB_2615052

Clone: 1MOD-1A9

Isotype: IgG1

Application(s): ICC, IF, IP, WB

Reactivity: Human, Mouse

Quantity: 100 μ g

Purification: Protein A Chromatography

Host: Mouse

Concentration: 0.60 μ g/ μ l

Molecular Weight: 25 kDa

Background: HP1 beta (Heterochromatin Protein 1 beta, Chromobox protein 1, CBX1) is one of the mammalian homologues of Drosophila HP1, first identified through its role in the maintenance of regions of the genome in a heterochromatic state. HP1 beta contains a chromo domain and a chromo shadow domain. Through the chromo domain, HP1 proteins are recruited to regions of chromatin that exhibit methylation of lysine 9 of histone H3. Interaction between HP1 beta and the lamin B receptor can facilitate the sequestration of heterochromatin to the nuclear envelope.

Immunogen: This HP1 β antibody was raised against recombinant mouse HP1 β .

Buffer: Purified IgG in 60 mM Sodium Citrate, 105 mM Tris-HCl, 30% glycerol, and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.